

# High Resolution

OEM weigh cells

WZA245-NC

WZA26-HC



WZA...	245-NC	26-HC
Capacity	240 g	20 g
Readability	10 µg	1 µg

The WZA-NC series is the successor of the well established WZA...CW-line. The products provide a readability of 1 µg | 10 µg for the weighing capacity of 20 g | 240 g. The weighing system as well the electronic are designed for space sensitive applications in stainless steel housings.

The weigh cell consist of two parts:

- A monolithic weighing system in stainless steel IP44 housing
- Stainless steel electronic box, all connection via the front

- Force compensation technology system with a resolution of 1 µg | 10 µg
- Superb electronics for 24 Mio. step resolution
- Internal calibration weight for easy checks in automated environments
- Load receptor with overload protection
- Possibility for under floor weighing
- Fast warm-up by separation of weighing system and electronics
- USB | RS232C interface for configuration and data transfer
- Optional PC-software with menu setting, weigh data display and internal calibration functions
- Applications can be found in weighing of samples in automated system, manual pipette calibration and in automated liquid control, e.g. LCD productions.

# Technical Data



Electronic box



YAC01CU (optional)



YCP01WZA (optional)

Model	WZA245-NC	WZA26-HC
Capacity	240 g	20 g
Readability	10 µg	1 µg
Preload <sup>1)</sup>	0 g	12.5 g
Repeatability (standard deviation) <sup>2)</sup>	≤± 20 µg	≤± 2 µg
Linearity	≤± 0.15 mg	≤± 0.02 mg
Measuring time <sup>2)3)</sup>	1.8 s   ± 10 µg	1.8 s   ± 1 µg
Operating temperature range	+10 ... +30 °C	
Allowable temperature range	+5 ... +40 °C	
Weigh cell dimension	153 × 232 × 75 mm (base × height)	
Electronic   PCB dimension	196 × 142 × 42 mm (base × height)	
Load receptor (with overload protection)	36 mm (diameter), weight of the load receptor is 13.5 ± 0.5 g	
Under floor load receptor	M3 thread	
Cable length weigh cell to electronics	2.8 m	
Power supply	TNG10   6971987; 115   230V <sub>AC</sub> , +15 ... -20%, 48 - 60Hz	
Power supply (alternative)	15...26, optimal 15V (+15% - 10%); 0.5V <sub>PP</sub> (Peak to Peak)	
Power consumption	9 VA (average)	3.75 VA (average)
Interface	USB-B, RS232C, software   hardware handshake; No system power via RS232 connector	

## Options | Accessories

Windows configuration software for test and adjustment	Sartorius CAS Suite
Connection cable 25   9 RS232   RS232-USB	RS232 7357314   YCC01-USBM2
Connection cable USB A to USB B	buy local
Liquid-crystal display	YAC01CU
Pipette calibration set	YCP01WZA

- 1) the preload can be configured by using the Sartorius CAS Suite software; higher preload will reduce the weighing range  
 2) depending on measurement equipment set-up and conditions;  
 3) measuring time is the time in which the measured value is in the given range around the static end value; test weight is approx. 25% of the weighing range.



**Germany**

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

**France & Suisse Romande**

Phone +33 170 62 50 00

**Finland & Baltics**

Phone +358 9 755 951

**Italy**

Phone +39 0362 5557 11

**U.K.**

Phone +44 1372 737159

**USA**

Toll-free +1 800 635 2906

**China**


Phone +86 21 6878 2300

**Japan**

Phone +81 3 6478 5200

**South Korea**

Phone +82 31 622 5700

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)